

PPACK

Thrombin inhibitor

Potent irreversible thrombin inhibitor. Inhibits thrombin activation of progelatinase A. Pretreatment with PPACK inhibits thrombin-induced cerebral thromboembolism *in vivo*.

Citations: 7

[View Online »](#)

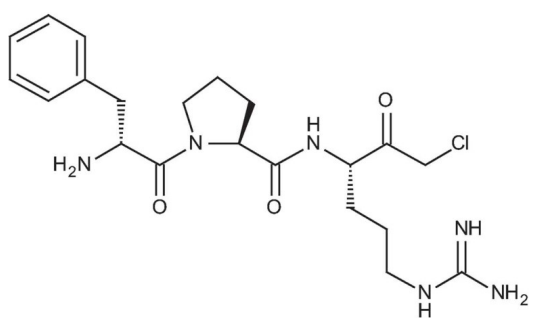
Ordering Information

[Order Online »](#)

BML-PI117-0005	5mg
BML-PI117-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	D-Phe-Pro-Arg-CMK
Appearance	White to off-white solid.
CAS	71142-71-7
Formula	$C_{21}H_{31}ClN_6O_3$
MW	451.0
Purity	≥90% (HPLC)
Sequence	D-Phe-Pro-Arg-chloromethylketone
Solubility	Soluble in water.

Technical Info / Product Notes Note: Product is not sterile.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)